

CERTIFICATE OF ANALYSIS

10X ELISA Wash Buffer (WB1)			
Catalog#	Size	Lot#	Expiry at room temperature
527	Bulk	16N52	06/2018
6251	25 mL	NA	NA
650	100 mL	NA	NA
651	500 mL	NA	NA
652	1 L	NA	NA
676	10 L	16N53	06/2018

Specifications	Criteria	Lot Data	Results
pH at 10X	6.8 – 7.0	6.89	Pass
pH at 1X	7.2 – 7.6	7.46	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	99,000 - 121,000 uMHO	115,833	Pass
Storage	Room temperature or refrigerated	Room temperature	Pass

Quality Control

Tracy Murphy
 Tracy Murphy

Date

06/23/2016
 06/23/2016